

5-Propargylamino- dCTP-XX-Cy5 (L pack)

Ordering Information

[Order Online »](#)

JBS-NU-809-XX-CY5-L

5x10 μ l

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance blue solution

Concentration 1.0 mM – 1.1 mM

Couple Type Modified nucleotides

Emission Maximum 670 nm

Excitation Maximum 649 nm

Formula $C_{57}H_{79}N_8O_{22}P_3S_2$ (free acid)

Formulation solution in 10 mM Tris-HCl

MW 1385.33 (free acid)

Purity $\geq 95\%$ (HPLC)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com